

PRODUCT INFORMATION



TL4-12

Item No. 42907

CAS Registry No.: 1620820-12-3
Formal Name: 4-methyl-3-[[6-(methylamino)-4-pyrimidinyl]oxy]-N-[3-(4-methyl-1-piperazinyl)-5-(trifluoromethyl)phenyl]-benzamide

MF: C₂₅H₂₇F₃N₆O₂

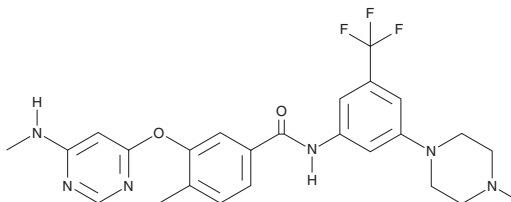
FW: 500.5

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

TL4-12 is supplied as a solid. A stock solution may be made by dissolving the TL4-12 in the solvent of choice, which should be purged with an inert gas. TL4-12 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

TL4-12 is an inhibitor of MAP4K2 (IC₅₀ = 37.2 nM).¹ It is selective for MAP4K2 over 20 other kinases at 1 μM. TL4-12 (500 nM) inhibits IL-1α- or TGF-β-induced increases in p38 phosphorylation in primary mouse embryonic fibroblasts. It induces apoptosis in MM.1S multiple myeloma cells when used at concentrations of 1 or 3 μM.²

References

1. Tan, L., Nomanbhoy, T., Gurbani, D., *et al.* Discovery of type II inhibitors of TGFβ-activated kinase 1 (TAK1) and mitogen-activated protein kinase kinase kinase 2 (MAP4K2). *J. Med. Chem.* **58**(1), 183-196 (2015).
2. Li, S., Fu, J., Yang, J., *et al.* Targeting the GCK pathway: A novel and selective therapeutic strategy against RAS-mutated multiple myeloma. *Blood* **137**(13), 1754-1764 (2021).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/27/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM